

TREVIGEN® Product Data

For Research Use Only. Not For Use In Diagnostic Procedures

Poly(ADP-ribose) (PAR) Polymer

Catalog #: 4336-100-01

Volume: 100 µl

Concentration: 10 µM

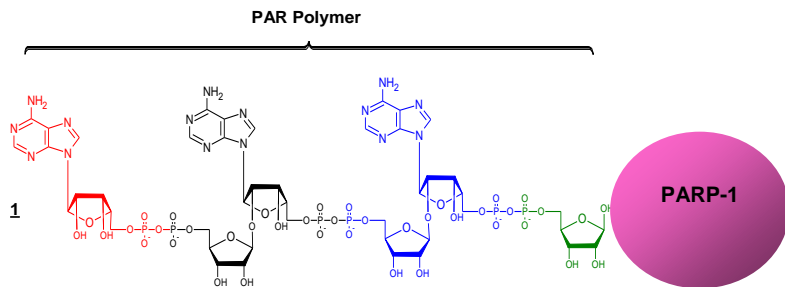


Figure 1: Poly (ADP-ribose) (PAR)

Description: PAR (figure 1) was synthesized using poly(ADP-ribose) polymerase (PARP-1) in the presence of NAD, cleaved from PARP-1, and purified by gel filtration. The PAR chain length ranges from 2 to 300 monomers, and it is recognized by Trevigen's PAR polymer monoclonal antibody (cat# 4335-MC-100) and polyclonal antibodies (cat# 4336-BPC-100). The PAR concentration was determined using the following equation:¹

$$[\text{PAR}] = \frac{(A_{258}) \text{ cm}^{-1}}{13,500 \text{ cm}^{-1} \text{ M}^{-1}}$$

Storage Conditions: PAR is provided in 10 mM Tris-HCl (pH 8.0), 1 mM EDTA and should be stored at -80 °C. It may be aliquoted to avoid repeated freeze-thawing.

Applications: Immunodetection as a positive control in ELISA and Western blot analysis.

References:

1. Shah, G.M., et al. 1995. Methods for biochemical study of poly(ADP-ribose) metabolism *in vitro* and *in vivo*. *Anal Biochem* **227**:1-13.
2. Affar, E.B., et al. 1998. Immunodot blot method for the de-tetection of poly(ADP-ribose) synthesized *in vitro* and *in vivo*. *Anal Biochem* **259**:280-283.
3. Menard, L. and G.G. Poirier. 1987. Rapid assay of poly (ADP-ribose) glycohydrolase. *Biochem Cell Biol* **65**:668-673.

TREVIGEN®

8405 Helgerman Court, Gaithersburg, MD 20877 USA

Voice: 1-800-TREVIGEN (1-800-873-8443) • 301-216-2800

Fax: 301-560-4973 • e-mail: info@trevigen.com • www.trevigen.com

**Poly(ADP-ribose)
(PAR) polymer**

Catalog#: 4336-100-01

Storage: -80 °C

TREVIGEN[®]
1-800-873-8443